

Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)

Non-reactive antibody to use as a negative control for mouse IgG1 primary antibodies. Available with CF® Dyes and other labels.

Product Description

This is a non-reactive mouse monoclonal IgG1 kappa antibody labeled with one of our bright and photostable CF® Dyes and other labels. This antibody may serve as a control for a mouse monoclonal IgG1 antibody under identical experimental conditions. Biotium also carries mouse monoclonal isotype control antibodies for [IgG2a Kappa \(MG2a-53\)](#) and [IgG2b Kappa \(IGG1/1333\)](#).

- Isotype control for antibody-based experiments
- Available with bright and photostable CF® Dyes
- APC, R-PE, RPE-Astral™616, APC-Astral™813, and biotin conjugates also available
- Suitable for western, flow cytometry, ELISA, immunofluorescence, and immunohistology

CF® Dyes offer exceptional brightness and photostability; learn more about [CF® Dyes](#).

Product attributes

Antibody number	#1331
Antibody reactivity (target)	Isotype Control
Antibody type	Isotype Control
Host species	Mouse
Clonality	Monoclonal
Clone	IGG1/1331
Isotype	IgG1, kappa
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@405S	411/431 nm	250 uL (25 ug) 1 mL (100 ug)	BNC041331-250 BNC041331-1ML	CF@405S Features
CF@405L	413/547 nm	250 uL (25 ug) 1 mL (100 ug)	BNC061331-250 BNC061331-1ML	CF@405L Features
CF@488A	562/584 nm	250 uL (25 ug) 1 mL (100 ug)	BNC881331-250 BNC881331-1ML	CF@488A Features
CF@503R	503/532 nm	250 uL (25 ug) 1 mL (100 ug)	BNC501331-250 BNC501331-1ML	CF@503R Features
CF@514	516/549 nm	250 uL (25 ug) 1 mL (100 ug)	BNC511331-250 BNC511331-1ML	CF@514 Features
CF@568	562/584 nm	250 uL (25 ug) 1 mL (100 ug)	BNC681331-250 BNC681331-1ML	CF@568 Features
CF@575	575/601 nm	250 uL (25 ug) 1 mL (100 ug)	BNC571331-250 BNC571331-1ML	
CF@594	593/615 nm	250 uL (25 ug) 1 mL (100 ug)	BNC941331-250 BNC941331-1ML	CF@594 Features
RPE-Astral™616	496, 546, 566/617 nm	250 uL (25 ug) 1 mL (100 ug)	BNC161331-250 BNC161331-1ML	
CF@640R	642/663 nm	250 uL (25 ug) 1 mL (100 ug)	BNC401331-250 BNC401331-1ML	CF@640R Features
CF@647	652/668 nm	250 uL (25 ug) 1 mL (100 ug)	BNC471331-250 BNC471331-1ML	CF@647 Features
CF@647Plus	652/668 nm	250 uL (25 ug) 1 mL (100 ug)	BNC641331-250 BNC641331-1ML	
CF@660C	667/685 nm	250 uL (25 ug) 1 mL (100 ug)	BNC601331-250 BNC601331-1ML	CF@660C Features
CF@700	696/721 nm	250 uL (25 ug) 1 mL (100 ug)	BNC001331-250 BNC001331-1ML	CF@700 Features
APC-Astral™813	633, 638/813 nm	250 uL (25 ug) 1 mL (100 ug)	BNC131331-250 BNC131331-1ML	
R-PE	496, 546, 565/576 nm	250 uL (25 ug) 1 mL (100 ug)	BNCR1331-250 BNCR1331-1ML	
APC	651/660 nm	250 uL (25 ug) 1 mL (100 ug)	BNCA1331-250 BNCA1331-1ML	
Biotin	N/A	250 uL (25 ug) 1 mL (100 ug)	BNCB1331-250 BNCB1331-1ML	
Purified	N/A	250 uL (50 ug)	BNUB1331-250	
Purified, BSA-free	N/A	50 uL (50 ug)	BNUM1331-50	

